

MULTISCREEN™ STABLE CELL LINE RAT RECOMBINANT BB1 RECEPTOR

PRODUCT INFORMATION

Catalog Number: C1211-1

Lot Number: C1211-1-121610

Quantity: 1 vial (2×10^6) frozen cells

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: C6

Transfection: Full-length rat NMBR cDNA (GenBank Accession Number NM_012799) with flag tag sequence at N-terminus

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM, 10% FBS

Stability: Stable after minimum of two months continuous growth

Data sheet

Background: The BB1 receptor (also known as NMB-R or Neuromedin B receptor) is a receptor for neuromedin-B (NMB), which is a mammalian bombesin-like peptide distributed widely in the central nervous system. The NMBR pathway is involved in the regulation of a wide variety of behaviors, such as spontaneous activity, feeding and anxiety-related behavior. A study using NMBR-deficient mice suggested that dysfunction in the NMBR pathway may constitute one of the risk factors of stress vulnerability.

Application: Functional assays

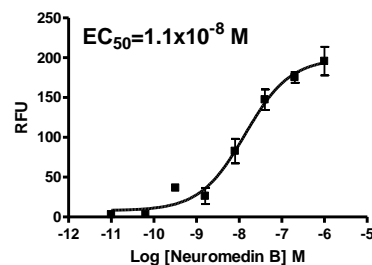


Figure legend: Dose-dependent calcium flux upon treatment with ligand, monitored with FlexStation.

References:

Moody *et al.* (2000) Nonpeptide neuromedin B receptor antagonists inhibit the proliferation of C6 cells. *Eur J Pharmacol* 409:133-142.

Yamada *et al.* (2002) Restraint stress impaired maternal behavior in female mice lacking the neuromedin B receptor (NMB-R) gene. *Neurosci Lett* 330:163-166.

Herold *et al.* (2003) The neuromedin B receptor antagonist, BIM-23127, is a potent antagonist at human and rat urotensin-II receptors. *Br J Pharmacol* 139:203-207.

FOR RESEARCH USE ONLY.

© 2005 Multispan Inc. All rights reserved. No part of this document may be reproduced in any form without prior permission in writing.

www.multispaninc.com
sales@multispaninc.com
support@multispaninc.com

Ver. October 2005

Phone: +1 (510) 887-0817
Fax: +1 (510) 887-0863
26219 Eden Landing Road
Hayward, CA 94545-3718
U.S.A.